

Product Number:	102732
Lot Number:	502333
Expiration:	24-Mar-2021
Platform(s):	HD-1/HD-X & SR-X



Data below represents results generated on the Simoa™ HD-1 Analyzer.

Component	Part Number	Lot Number
Bead Reagent	102722	010508
Detector Reagent	102724	010404
SBG Reagent	102726	010403
Sample Diluent	102728	010407
Calibrator Diluent	102731	010406
Calibrator Concentrate	100658	920712
RGP Reagent	103159	N/A <sup>1</sup>

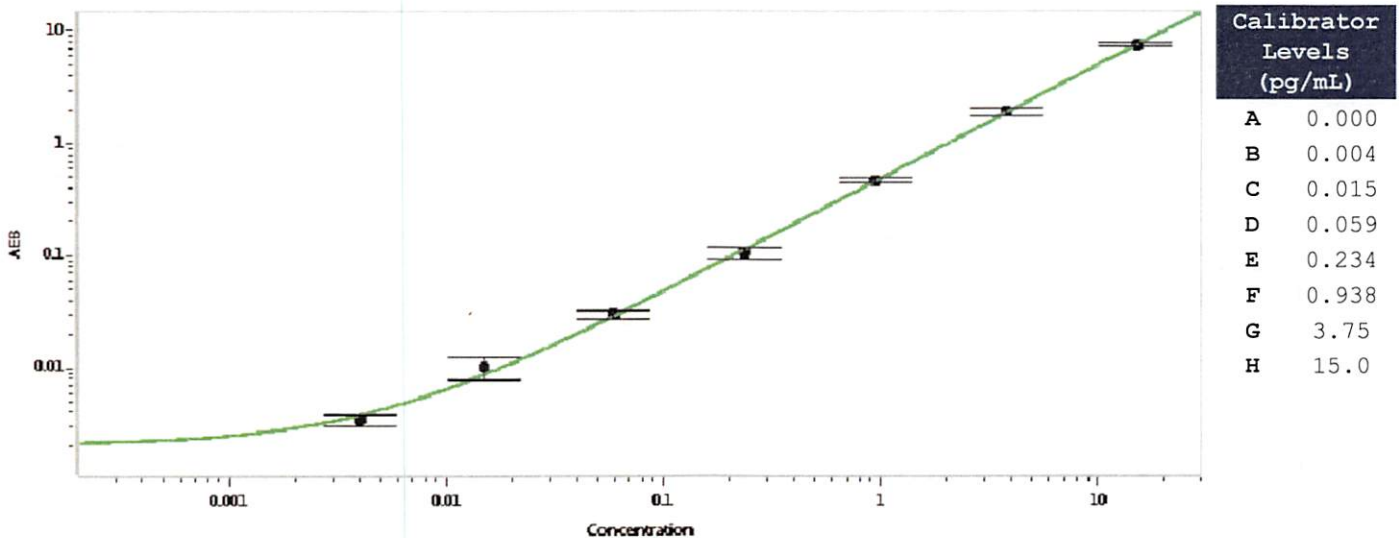
<sup>1</sup> RGP is not Kit Lot Specific

Release Materials*			
	Lot	Result (pg/mL)	Mean Range (pg/mL)
Control 1	909502	0.155	0.095-0.162
Control 2	909502	3.78	2.88-4.52
Panel 1	911401	0.032	0.013-0.050
Panel 3	911403	1.76	1.22-2.02
Panel 4	911404	5.69	3.95-7.43

\* Ranges shown are generated internally for new lot release only. Customer should generate their own control ranges.

Calibrator Concentrate Value Assignment (pg/mL)	412
---	-----

Calibration Curve



Review/Approval	<u>Shivani Goel</u> QC Manager	<u>Shivani Goel</u> 16 Apr 2020
	Name	Signature/Date